

**CERTIFICATE OF ANALYSIS**

Control Pack Description: Chemokine Panel 1 (human) Control Pack
Catalog Number: C4047-1
Lot Number: CP000194
Expiration Date: 31 JAN 2031

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Control 1	A00C1061	2-8°C	31 MAY 2031
Chemokine Panel 1 (human) Control 2	A00C1062	2-8°C	31 MAY 2031
Chemokine Panel 1 (human) Control 3	A00C1047	2-8°C	31 JAN 2031

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
Eotaxin	1730	596	182	pg/mL
MIP-1 β	735	119	22.9	pg/mL
Eotaxin-3	4200	638	79.7	pg/mL
TARC/CCL17	1460	231	34.5	pg/mL
IP-10	806	120	17.6	pg/mL
MIP-1 α	1040	440	177	pg/mL
IL-8 (HA)	66200	22900	8230	pg/mL
MCP-1	562	83.9	13.5	pg/mL
MDC	4250	1840	857	pg/mL
MCP-4	666	212	61.7	pg/mL

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		01 SEP 2026